



Certificate of Analysis

CloneDetect

Human IgG (H+L) Specific, Fluorescein

BSA Free

K8202

Batch #	CDT0317A
----------------	-----------------

Full Description	Sheep anti-human IgG (H+L), polyclonal, Fluorescein-conjugated. Sterile, BSA and azide free
Origin of Host Species	USA Sheep kept at a registered technical plant (1774/2002 registration # VTTEC-0002) under regular veterinary supervision.
Immunogen	Human IgG (H+L) purified from human serum
Preparation	Antiserum is prepared by immunising sheep with the above immunogen. The antiserum is then conjugated with fluorescein. Unreacted fluorochrome is removed by gel filtration.
Format	Fluorescein-conjugated IgG, liquid
Specificity	The unconjugated immunoglobulin was shown to be reactive to human serum and unreactive to mouse/rate/bovine serum by immunoelectrophoresis
Method of Purification	Affinity purification by absorption to immobilised human IgG
Viral Testing	Negative. Product was found to be free of viral contaminants based on: qualified real-time PCR screening for rodent virus panels. Tests performed by Charles River (full report available on request).
Mycoplasma Testing	Negative. Product is assayed following a broth enrichment step using an RT-PCR assay. Assay meets or exceeds the requirements of 21 CFR 610.30, 1993 PTC, and EP. Test performed by Charles River (report available on request).
Appearance	Clear yellow solution
Concentration	10,000 U/ml
Volume	1.0 ml
Total Amount	0.75 mg
Dye : Protein Ratio	> 4.5
Buffer	Phosphate Buffered Saline pH7.2

Sterility	0.2µm sterile filtered
Expiry Date	12 months from receipt of product
Storage	2 – 8°C
Preservatives/Stabilizers	No BSA, No Azide

This product has been manufactured to ISO9001:2008 Quality Assurance guidelines

For research use only

Signed: *Kieu-Khiem Nguyen*

Date: 10 – MAY - 2018

Production Biologist